

Riluzole

Na⁺ channel blocker, K⁺ channel activator

Glutamate release inhibitor. Neuroprotective. Blocks the inactivated state of voltage-gated neuronal Na⁺ channels at very low concentrations (about 0.1µM) and activates the TREK family of two-pore domain potassium channels.

Citations: 11

[View Online »](#)

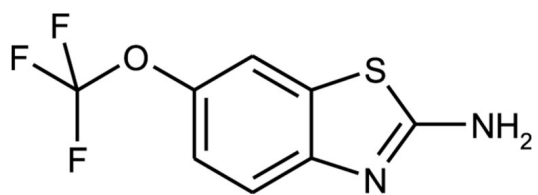
Ordering Information

[Order Online »](#)

BML-NA138-0020	20mg
----------------	------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Stable for up to 1 year after receipt when stored at +20°C. Stock solutions are stable for up to 3 months at -20°C.
Long Term Storage	Ambient
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	2-Amino-6-(trifluoromethoxy)-benzothiazole, 6-(Trifluoromethoxy)-2-benzothiazolamine
Appearance	White to off-white solid.
CAS	1744-22-5
Couple Target	Potassium channel, Sodium channel
Couple Type	Activator, Blocker
Formula	$C_8H_5F_3N_2OS$
MI	14: 8223
MW	234.2
Purity	≥98% (TLC)
Solubility	Soluble in DMSO (>25mg/ml), 100% ethanol (>25mg/ml), methanol (>40mg/ml), 45% (w/v) aq 2-hydroxypropyl-β-cyclodextrin (14mg/ml) or in 0.1N HCl (11mg/ml); insoluble in water.

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-fr@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com